

PRODUCT DATASHEET
ChemiScreen™ CCR1 Chemokine Membrane Preparation

CATALOG NUMBER: HTS005M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

BACKGROUND: CCR1 is a GPCR that binds to a variety of CC ligands, including MIP-1 α , RANTES, MCP-3, HCC-1, HCC-2, HCC-4, and MPIF-1 (Olson and Ley, 2002). Lymphocytes, macrophages, dendritic cells, and GM-CSF-activated neutrophils express CCR1 (Kaufmann *et al.*, 2001; Cheng *et al.*, 2001). Two selective, non-peptide small molecule antagonists of CCR1, BX-471 and CP-481,715 have been synthesized (Gladue *et al.*, 2003; Liang *et al.*, 2000). Pharmacological and genetic targeting of CCR1, either alone or in combination with cyclosporin A, reduces cardiac and renal allograft rejection (Gao *et al.*, 2000; Horuk *et al.*, 2001a; Horuk *et al.*, 2001b), allergic encephalomyelitis (Liang *et al.*, 2000), and renal fibrosis (Anders *et al.*, 2002) in experimental models. CCR1 Membrane Preparations are ideal for screening molecules that disrupt interactions between CCR1 and its ligands. The CCR1 Membrane Preps bind to MIP-1 α with a K_d of 1.8 nM, and yield greater than a 10-fold signal-to-background ratio with [¹²⁵I]-MIP-1 α at a concentration of 0.1 nM.

APPLICATIONS: Radioligand Binding Assay

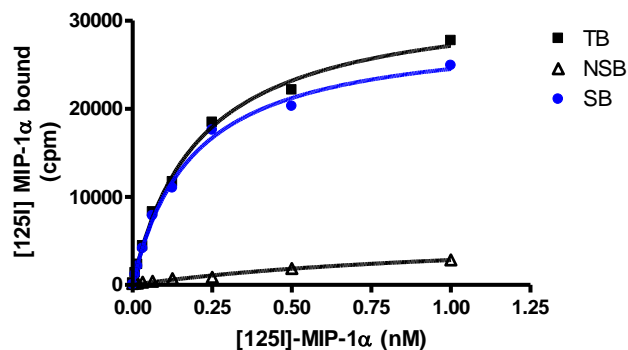


Figure 1. Saturation Binding for CCR1. 5 μ g/well of CCR1 Membrane Preparation were incubated with increasing amounts of [¹²⁵I]-MIP-1 α in the absence (total binding, TB) or presence (nonspecific binding, NSB) of an excess of unlabeled MIP-1 α . Specific binding (SB) was determined by subtracting NSB from TB. The sample data are from a representative lot.

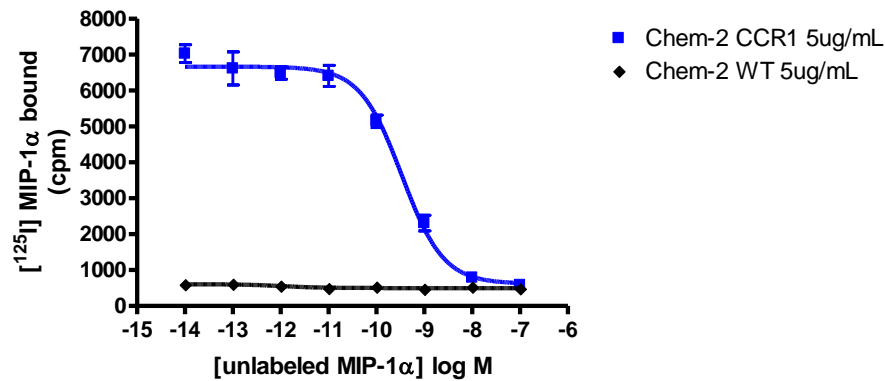


Figure 2. Competition binding of MIP-1 α to CCR1 Membrane Preparations. CCR1 Membrane Preparation (5 μ g/well) or Wild-Type Chem-2 Membrane Preparation (catalog # HTS000MC2) at 10 μ g/well were incubated with 0.1 nM [125 I]-MIP-1 α and increasing concentrations of unlabeled MIP-1 α . Greater than 10-fold signal to background was obtained with the CCR1 Membrane Preparation, whereas negligible signal was obtained with the Wild-Type Chem-2 Membrane Preparation. Representative sample data.

SPECIFICATIONS: 1 unit = 5 μ g
 B_{max} for [125 I]MIP-1 α binding: 3.3 pmol/mg protein
 K_d for [125 I] MIP-1 α binding: ~ 0.02 nM
Signal:background: >10-fold

TRANSFECTION: Full-length human CCR1 cDNA (Accession Number: L09230)

HOST CELLS: Chem-2.

RECOMMENDED ASSAY CONDITIONS: Membranes are mixed with radioactive ligand and unlabeled competitor (see Figures 1 and 2 for concentrations tested) in binding buffer in a nonbinding 96-well plate, and incubated for 1-2 h. Prior to filtration, a GF/C 96-well filter plate is coated with 0.33% polyethyleneimine for 30 min, then washed with 50mM HEPES, pH 7.4, 0.5% BSA. Binding reaction is transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The plate is dried and counted.

Binding buffer: 50 mM Hepes, pH 7.4, 5 mM MgCl₂, 1 mM CaCl₂, 0.2% BSA, filtered and stored at 4°C

Radioligand: [125 I] MIP-1 α (Perkin Elmer # NEX298)

Wash Buffer: 50 mM Hepes, pH 7.4, 500mM NaCl, 0.1% BSA, filtered and stored at 4°C.

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 10-fold signal:background with 125 I-labeled MIP-1 α at 0.1 nM.

PRESENTATION: Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives.
Packaging method: Membrane proteins were adjusted to the indicated concentration in 1 ml packaging buffer, rapidly frozen, and stored at -80°C.

STORAGE/HANDLING: Store at -70°C. Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

REFERENCES:

1. Anders, H.J., *et al.* (2002) A chemokine receptor CCR-1 antagonist reduces renal fibrosis after unilateral ureter ligation. *J. Clin. Invest.* 109: 251-9.
2. Cheng, S.S., *et al.* (2001) Granulocyte-macrophage colony stimulation factor up-regulates CCR1 in human neutrophils. *J. Immunol.* 166: 1178-84.
3. Gao, W., *et al.* (2000) Targeting of the chemokine receptor CCR1 suppresses development of acute and chronic cardiac allograft rejection. *J. Clin. Invest.* 105: 35-44.
4. Gladue, R.P., *et al.* (2003) CP-481,715, a potent and selective CCR1 antagonist with potential therapeutic implications for inflammatory diseases. *J. Biol. Chem.* 278: 40473-80.
5. Horuk, R., *et al.* (2001a) CCR1-specific non-peptide antagonist: efficacy in a rabbit allograft rejection model. *Immunol. Lett.* 76: 193-201.
6. Horuk, R., *et al.* (2001b) A non-peptide functional antagonist of the CCR1 chemokine receptor is effective in rat heart transplant rejection. *J. Biol. Chem.* 276: 4199-4204.
7. Kaufmann, A., *et al.* (2001) Increase of CCR1 and CCR5 expression and enhanced functional response to MIP-1 alpha during differentiation of human monocytes to macrophages. *J. Leukoc. Biol.* 69: 248-52.
8. Liang, M., *et al.* (2000) Identification and characterization of a potent, selective, and orally active antagonist of the CC chemokine receptor-1. *J. Biol. Chem.* 275: 19000-8.
9. Olson, T.S. and Ley, K. (2002) Chemokines and chemokine receptors in leukocyte trafficking. *Am. J. Physiol. Regul. Integr. Comp. Physiol.* 283: R7-R28.

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services